

Product Name:	Exendin 4	
Catalog Number:	AS-24463 (0.5 mg) AS-24464 (1 mg)	Lot Number: See label on vial
Sequence:	H-His-Gly-Glu-Gly-Thr-Phe-Thr-Ser-Asp-Leu-Ser-Lys-Gln-Met-Glu-Glu- Glu-Ala-Val-Arg-Leu-Phe-Ile-Glu-Trp-Leu-Lys-Asn-Gly-Gly-Pro-Ser-Ser- Gly-Ala-Pro-Pro-Pro-Ser-NH2 (3-letter code) HGEGTFTSDLSKQMEEEAVRLFIEWLKNGGPSSGAPPPS-NH2 (1-letter code)	
Molecular Weight:	4187.6	
% Peak Area by HPLC: ≥ 95		
Appearance:	Lyophilized white powder	
Peptide Reconstitution: Exendin 4 peptide is freely soluble in H <sub>2</sub> O.		

Storage: Exendin 4 peptide is shipped at ambient temperature. Upon receipt, store lyophilized

peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20°C or lower.

Description: This peptide, like Exendin-3, stimulates an increase in acinar cAMP, without stimulating the release of amylase. Ref: Eng, J. et al. *J. Biol. Chem.* **267**, 7402 (1992).

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

Animals were followed for 8 days. They were divided into five groups: (1) 30 islets with exendin-4 treatment (AnaSpec; San Jose, CA, USA); (2) 30 islets without treatment; (3) 75 islets without treatment (quality controls); (4) non-transplanted diabetic mice treated with exendin-4; and (5) non-transplanted diabetic mice without treatment. Exendin-4 was given as a single i.p. injection of 100 ng (approximately 24 pmol) per day from day 0 to day 7-<u>Sharma, A. et al.</u> <u>Diabetologia 49, 1247 (2006).</u>

Administration of exendin-4 Synthetic exendin-4 (Ex-4) (AnaSpec, USA) was dissolved in PBS. For the glucose tolerance and insulin secretion tests, Ex-4 was administered i.p. at a dose of 0.1g/g body weight 30 minutes-<u>Ding,J. et al. *Endocrinology* 10.1210/en.2008-1006 (2009).</u>

**Published Citations:** 

<u>Sharma, A. et al. *Diabetologia* **49**, 1247 (2006).</u> <u>Ding,J. et al. *Endocrinology* 10.1210/en.2008-1006 (2009).</u>

For Research Use Only